

# 15-Deoxy- $\Delta$ 12,14-prostaglandin J2

## PPAR $\gamma$ ligand

Prostaglandin D2 derived nuclear peroxisome proliferator-activated receptor  $\gamma$  (PPAR $\gamma$ ) ligand. Induces apoptosis in a number of cell types including human lung cancer cells, synoviocytes and macrophages. Promotes differentiation of fibroblasts to adipocytes. Antiinflammatory activity. Inhibits angiogenesis. Prevents TNF $\alpha$  induced leptin increase in mice. TRPA1 agonist. Serious doubts as to the biological relevance of this prostaglandin have been raised. Endogenous levels seem to be too low to have any effect as a PPAR $\gamma$  ligand.

Citations: 18

[View Online »](#)

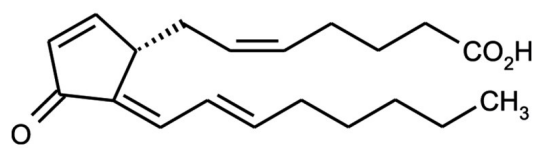
## Ordering Information

[Order Online »](#)

BML-PG050-0001	1mg
----------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
<b>Handling</b>	It is recommended to add an appropriate solvent directly to the vial.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	15-deoxy- $\Delta$ 12,14-PGJ2
<b>Appearance</b>	Neat oil. Vials may appear empty because the oil tends to coat the inside of the vial.
<b>CAS</b>	87893-55-8
<b>Couple Target</b>	PPAR, TRP channel
<b>Couple Type</b>	Ligand
<b>Formula</b>	$C_{20}H_{28}O_3$
<b>MW</b>	316.5
<b>Purity</b>	$\geq$ 97% (TLC)
<b>Solubility</b>	Soluble in 100% ethanol (25mg/ml) or DMSO (25mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)